
CYP3A7 Polyclonal Antibody, FITC Conjugated

(Catalog A62368)

Background

Cytochromes P450 are a group of heme-thiolate monooxygenases. In liver microsomes, this enzyme is involved in an NADPH-dependent electron transport pathway. It oxidizes a variety of structurally unrelated compounds, including steroids, fatty acids, and xenobiotics.

Description

CYP3A7 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P24462

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cytochrome P450 3A7 protein (1-200AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Cytochrome P450 3A7, CYP3A7, Cytochrome P450-HFLA, CYP3A7

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
